

# IL-6 High Sensitivity Controls

## Human IL-6 controls for use with the IL-6 (human) high sensitivity ELISA kit

This product is a set of quantitative controls for the determination of IL-6 concentration in biological matrices, such as serum and plasma. The concentrations have been determined using the Enzo Life Sciences IL-6 (human) high sensitivity ELISA kit (ENZ-KIT178). The controls were lyophilized in porcine serum and contain human IL-6 at a low, medium and high concentration.

Interleukin-6 (IL-6) is a pleiotropic pro-inflammatory cytokine, involved in the induction, growth, and differentiation of cells in the immune and hematopoietic systems, and initiates and coordinates inflammatory reaction. Formerly known as B-cell differentiation factor, IL-6 is a 24 kDa protein that is 212 amino acids in length. It is produced by various types of cells, such as macrophages, T-cells, B-cells and fibroblasts. IL-6 is one of the essential factors for antibody production in B cells, a potent growth factor for myeloma and plasmocytoma cells, and induces acute phase proteins in hepatocytes and promotes expansion of activated T-cells. Levels of IL-6 have been shown to increase in serum as a result of burn trauma and sepsis and also in cerebrospinal fluid with acute viral or bacterial infections of the central nervous system (CNS). Increased levels of IL-6 are observed in several inflammatory diseases, such as rheumatoid arthritis, Castleman's disease and systemic juvenile idiopathic arthritis as well as playing a critical role in the pathogenesis of autoimmune disease. In addition to all of the functions described above, IL-6 also acts as a myokine, which is a cytokine produced by working muscle during exercise.

### Ordering Information

[Order Online »](#)

ENZ-KIT178-CTL	1Set
----------------	------

### Manuals, SDS & CofA

[View Online »](#)

- Lyophilized set of controls
- Includes high, medium, and low controls

## Handling & Storage

**Short Term Storage** -20°C

**Long Term Storage** -80°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Interleukin-6

**Application** ELISA

**Contents** 1 high, 1 medium, and 1 low control

**Formulation** Lyophilized in porcine serum.

**Reconstitution** Each control should be reconstituted in 0.5 mL deionized or distilled water.

**UniProt ID** P05231



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)